

4-Phenylchalcone Oxide

Epoxide hydrolase inhibitor

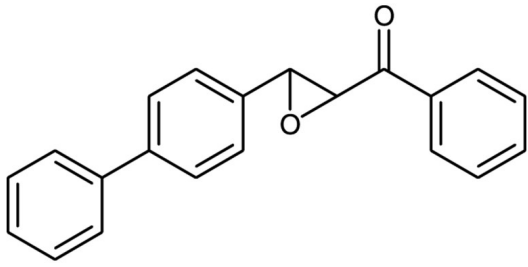
Potent and selective inhibitor of microsomal and cytosolic epoxide hydrolase (IC_{50} =68 nM). Inhibits epoxide hydrolysis of epoxyeicosatrienoic acids.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to one year. Store solutions at -20°C for up to three months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	82389-34-2
Couple Target	Epoxide hydrolase
Couple Type	Inhibitor
Formula	$C_{21}H_{16}O_2$
MW	300.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml)
Source	Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com